

PRODUCT INFORMATION



ANGPTL3 (human, recombinant)

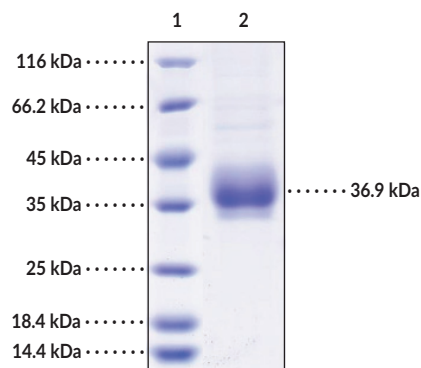
Item No. 42772

Overview and Properties

Synonyms: ANG-5, Angiopoietin-5, Angiopoietin-like 3, Angiopoietin-related Protein 3
Source: Recombinant human C-terminal His-tagged ANGPTL3 expressed in HEK293 cells
Amino Acids: 17-220
Uniprot No.: Q9Y5C1
Molecular Weight: 25.3 kDa
Storage: -80°C (as supplied)
Stability: ≥1 year
Purity: ≥95% estimated by SDS-PAGE
Supplied in: Lyophilized from sterile PBS, pH 7.4.
Endotoxin Testing: <1.0 EU/μg, determined by the LAL endotoxin assay

Information represents the product specifications. Batch specific analytical results are provided on each certificate of analysis.

Image



Lane 1: MW Markers
Lane 2: ANGPTL3

SDS-PAGE Analysis of ANGPTL3. This protein has a calculated molecular weight of 25.3 kDa. It has an apparent molecular weight of approximately 36.9 kDa by SDS-PAGE under reducing conditions due to glycosylation.

WARNING
THIS PRODUCT IS FOR RESEARCH ONLY - NOT FOR HUMAN OR VETERINARY DIAGNOSTIC OR THERAPEUTIC USE.

SAFETY DATA
This material should be considered hazardous until further information becomes available. Do not ingest, inhale, get in eyes, on skin, or on clothing. Wash thoroughly after handling. Before use, the user must review the [complete](#) Safety Data Sheet, which has been sent via email to your institution.

WARRANTY AND LIMITATION OF REMEDY
Buyer agrees to purchase the material subject to Cayman's Terms and Conditions. Complete Terms and Conditions including Warranty and Limitation of Liability information can be found on our website.

Copyright Cayman Chemical Company, 04/17/2025

CAYMAN CHEMICAL

1180 EAST ELLSWORTH RD
ANN ARBOR, MI 48108 · USA

PHONE: [800] 364-9897

[734] 971-3335

FAX: [734] 971-3640

CUSTSERV@CAYMANCHEM.COM

WWW.CAYMANCHEM.COM

PRODUCT INFORMATION



Description

Angiopoietin-like 3 (ANGPTL3) is an angiopoietin-related protein involved in lipid metabolism.^{1,2} It is composed of a signal peptide, coiled-coil domain, linker region, and C-terminal fibrinogen-like domain.^{2,3} ANGPTL3 is expressed in adult liver and during embryonic development.³ It inhibits the activity of lipoprotein lipase and endothelial lipase, which increases plasma triglycerides and HDL-cholesterol (HDL-C) levels.⁴⁻⁶ Loss-of-function mutations in ANGPTL3 are associated with decreased risk of coronary artery disease, coronary atherosclerosis, or myocardial infarction.⁶ Formulations containing monoclonal antibodies targeting ANGPTL3 have been used in the treatment of homozygous familial hypercholesterolemia. Cayman's ANGPTL3 (human, recombinant) protein consists of 215 amino acids and has a calculated molecular weight of 25.3 kDa. By SDS-PAGE, under reducing conditions, the apparent molecular mass of the protein is 36.9 kDa due to glycosylation.

References

1. Kadamatsu, T., Tabata, M., and Oike, Y. Angiopoietin-like proteins: Emerging targets for treatment of obesity and related metabolic diseases. *FEBS J.* **278**(4), 559-564 (2011).
2. Oike, Y., Akao, M., Kubota, Y., et al. Angiopoietin-like proteins: Potential new targets for metabolic syndrome therapy. *Trends Mol. Med.* **11**(10), 473-479 (2005).
3. Conklin, D., Gilbertson, D., Taft, D.W., et al. Identification of a mammalian angiopoietin-related protein expressed specifically in liver. *Genomics* **62**(3), 477-482 (1999).
4. Shan, L., Yu, X.C., Liu, Z., et al. The angiopoietin-like proteins ANGPTL3 and ANGPTL4 inhibit lipoprotein lipase activity through distinct mechanisms. *J. Biol. Chem.* **284**(3), 1419-1424 (2009).
5. Shimamura, M., Matsuda, M., Yasumo, H., et al. Angiopoietin-like protein3 regulates plasma HDL cholesterol through suppression of endothelial lipase. *Arterioscler. Thromb. Vasc. Biol.* **27**(2), 366-372 (2007).
6. Luo, F., Das, A., Khetarpal, S.A., et al. ANGPTL3 inhibition, dyslipidemia, and cardiovascular diseases. *Trends Cardiovasc. Med.* **34**(4), 215-222 (2024).

CAYMAN CHEMICAL
1180 EAST ELLSWORTH RD
ANN ARBOR, MI 48108 · USA
PHONE: [800] 364-9897
[734] 971-3335
FAX: [734] 971-3640
CUSTSERV@CAYMANCHEM.COM
WWW.CAYMANCHEM.COM